



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
ACE2 Inhibitor Screening Assay Kit
Catalog #79923
Size: 96 reactions

BACKGROUND: Angiotensin converting enzyme 2 (ACE2) is an exopeptidase that catalyzes the conversion of angiotensin II to angiotensin 1-7 and L-phenylalanine. Angiotensin II is part of the classical renin angiotensin system (RAS), a hormone system that regulates fluid balance, blood pressure and maintains vascular tone. ACE2 has been also proved to be the receptor for the human respiratory coronavirus NL63, the SARS-coronavirus (SARS-CoV) and the novel coronavirus 2019-nCoV/SARS-CoV-2.

DESCRIPTION: The *ACE2 Inhibitor Screening Assay Kit* is designed to measure the exopeptidase activity of ACE2 for screening and profiling applications. The ACE2 assay kit comes in a convenient 96-well format, with purified ACE2, its substrate, and ACE2 buffer for 96 reactions.

COMPONENTS:

Catalog #	Component	Amount	Storage	
11003	ACE2	2 µg	-80 °C	Avoid multiple freeze/thaw cycles!
	ACE2 Fluorogenic Substrate	3 ml	-20 °C	
	ACE2 Buffer	3 ml	-20 °C	
79685	96-well black microplate	1	Room Temp	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Adjustable micropipettor and sterile tips
Fluorescent microplate reader

APPLICATIONS: Useful for studying enzyme kinetics and screening small molecule inhibitors and antibodies for drug discovery and HTS applications.

STABILITY: Up to 6 months from date of receipt, when stored as recommended.

REFERENCES:

Danilczyk, U., *et al. Circulation Research*. 2006. **98(4)**: 463-471.
Ge, X. Y., *et al. Nature*. 2013. **503(7477)**: 535-538.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

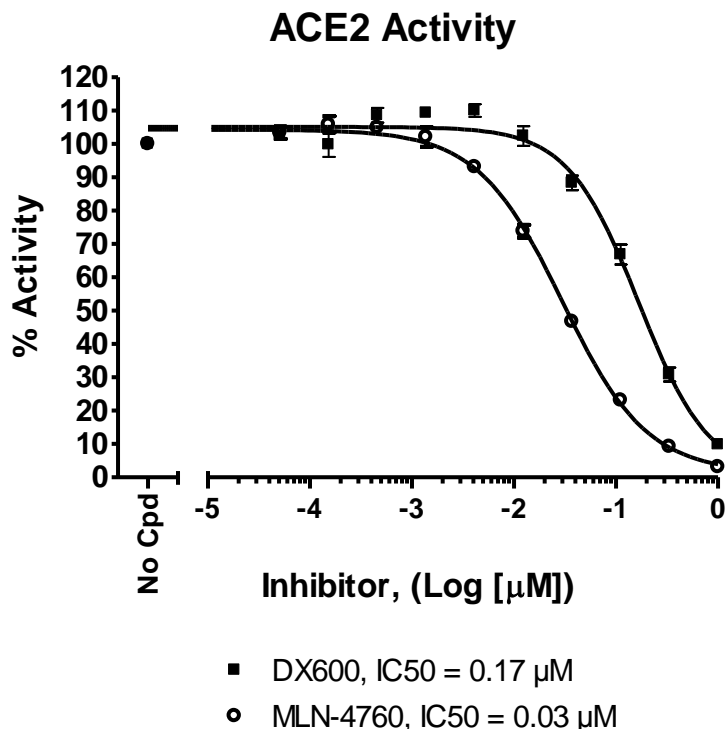
1. Thaw **ACE2** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Note: **ACE2** is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
2. Prepare **Enzyme solution** (0.5 ng/ μ l ACE2) by diluting ACE2 in **ACE2 Buffer**.
3. Add 20 μ l of **Enzyme solution** (0.5 ng/ μ l ACE2) to each well designated "Positive Control", and "Test Inhibitor" and 20 μ l of **ACE2 buffer** to each well designated "Blank".
4. Add 5 μ l of **Test Inhibitor** solution to each well designated "Test Inhibitor". For the wells labeled "Positive Control" and "Blank", add 5 μ l of **Inhibitor buffer** (same solution without inhibitor). Note: Keep the final concentration of DMSO in the reaction \leq 1%.
5. Add 25 μ l of **ACE2 Fluorogenic Substrate** to all wells. Protect from light and incubate reaction at room temperature for 60 minutes.

	Positive Control	Test Inhibitor	Blank
Enzyme solution (0.5 ng/ μ l ACE2)	20 μ l	20 μ l	-
ACE2 Buffer	-	-	20 μ l
Test inhibitor	-	5 μ l	-
Inhibitor buffer (no inhibitor)	5 μ l	-	5 μ l
ACE2 Fluorogenic Substrate	25 μ l	25 μ l	25 μ l
Total	50 μl	50 μl	50 μl

6. Read fluorescence intensity of the samples ($\lambda_{\text{excitation}} = 555 \text{ nm}$; $\lambda_{\text{emission}} = 585 \text{ nm}$) in an appropriate microplate reader.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com

Example of assay results:



ACE2 inhibition by DX600 (Cayman, #22186) and MLN-4760 (Sigma, #5306160001) measured using the *ACE2 Inhibitor Screening Assay Kit*, BPS Bioscience Catalog #79923. Fluorescence was measured using a Bio-Tek microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product</u>	<u>Catalog#</u>	<u>Size</u>
ACE2	11003	10 μ g
PrCP, His-tag	80380	10 μ g
Renin, Human, His-tag	80200	100 μ g
Renin, Rat, His-tag	80201	100 μ g
Renin Assay Kit	80211	96 rxns.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com